

# CHIMERY'S: An AI-Driven Leap Forward in Peptide Identification

Martin Frejno<sup>1</sup>; Daniel P Zolg<sup>1</sup>; Tobias Schmidt<sup>1</sup>; Siegfried Gessulat<sup>1</sup>; Michael Graber<sup>1</sup>; Florian Seefried<sup>1</sup>; Magnús Rathke-Kuhnert<sup>1</sup>; Samia Ben Fredj<sup>1</sup>; Shyamnath Premnadh<sup>1</sup>; Patroklos Samaras<sup>1</sup>; Kai Fritzsche<sup>2</sup>; Frank Berg<sup>2</sup>; Waqas Nasir<sup>2</sup>; David Horn<sup>3</sup>; Bernard Delanghe<sup>2</sup>; Christoph Henrich<sup>2</sup>; Bernhard Kuster<sup>4</sup>; Mathias Wilhelm<sup>4</sup> <sup>1</sup>MSAID GmbH, Garching b.München, Germany; <sup>2</sup>Thermo Fisher Scientific (Bremen) GmbH, Bremen, Germany; <sup>3</sup>Thermo Fisher Scientific, San Jose, CA; <sup>4</sup>Technical University of Munich, Freising, Germany

## ABSTRACT

**Purpose:** Chimeric spectra represent a substantial challenge for bottom-up proteomics data analysis. Here, we describe CHIMERY'S™, a novel, highly scalable, cloud-native, microservice-based and artificial intelligence-powered search algorithm that rethinks the analysis of tandem mass spectra from the ground up and deconvolutes chimeric spectra based on predicted fragment ion intensities.

**Methods:** We performed comparative analyses of standard HeLa tryptic digests that were acquired on various mass spectrometry platforms using different gradient lengths and isolation widths, as well as *in-silico* generated and publicly available datasets from various organisms using Sequest HT™, the Precursor Detector Node, INFERY'S™ Rescoring [1] and CHIMERY'S™ as implemented in a pre-release version of Thermo Scientific™ Proteome Discoverer™ 3.0 software.

**Results:** CHIMERY'S doubles peptide identifications in classical data-dependent acquisition (DDA) datasets compared to Sequest HT and increases the number of identified peptides per protein by 2.5-fold on average, which translates to ~2 PSMs per spectrum and an identification rate of >80%. Entrapment analyses suggest that the CHIMERY'S score set is well-calibrated and dilution experiments confirm that peptides unique to CHIMERY'S follow the expected ratio distribution. Experiments based on simulated chimeric spectra establish that CHIMERY'S has a sensitivity of >90%. Using CHIMERY'S enables more efficient data acquisition strategies, as both wider isolation windows and shorter gradients can be used to generate more PSMs in a shorter timeframe.

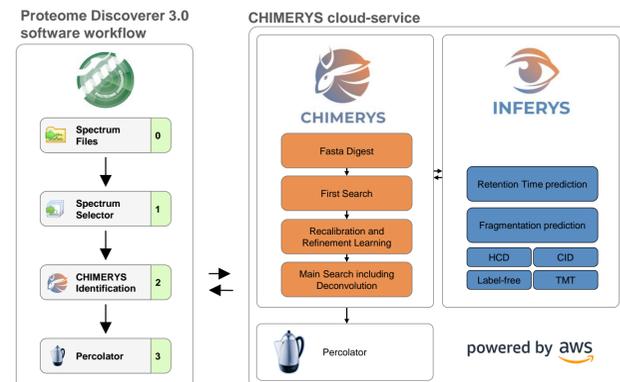
## INTRODUCTION

Matching peptide sequences to tandem mass spectra is integral to bottom-up proteomics. Chimeric spectra are estimated to constitute >40% of DDA data [2], violating the assumption that one spectrum represents one peptide. Some search engines allow multi-pass searches or duplicate chimeric spectra for several possible precursors, but few account for the fact that the measured intensities of (isobaric) fragment ions may be the sum of multiple peptides. This introduces errors and leaves valuable information unused, resulting in far fewer peptide identifications than contained in the data. Here, we describe CHIMERY'S, a new AI-based search algorithm that rethinks the analysis of tandem mass spectra from the ground up. It routinely doubles the number of peptide identifications in comparison to classical search algorithms and reaches identification rates of >80%.

## MATERIALS AND METHODS

### Data Analysis

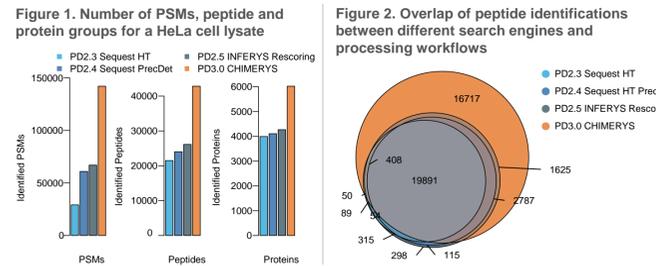
CHIMERY'S is a cloud-native search algorithm that uses accurate predictions of peptide fragment ion intensities and retention times provided by the deep learning framework INFERY'S 2.0. Based on an initial coarse search, INFERY'S performs data-driven model refinement to maximize prediction accuracy. Tandem mass spectra are analyzed without pre-processing or candidate selection using features detected in precursor mass spectra. Instead, all candidates in the isolation window of a given tandem mass spectrum are considered simultaneously and compete for measured fragment ion intensity in one concerted step. CHIMERY'S aims to explain as much measured intensity with as few candidate peptides as possible, resulting in the deconvolution of chimeric spectra. Peptide-spectrum match (PSM)-level false discovery rate (FDR)-control is performed using Percolator [3]. CHIMERY'S profits from cloud-based parallelization and is available through a node in a pre-release version Thermo Scientific Proteome Discoverer 3.0 software.



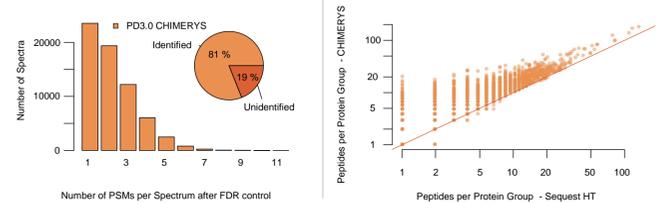
## RESULTS

### CHIMERY'S doubles peptide identifications in single-shot full proteome DDA datasets

CHIMERY'S deconvolution algorithm identifies peptides hidden in chimeric spectra of DDA data files. Here, a digest of a HeLa cell lysate was analyzed using a 1-hour gradient on a Thermo Scientific™ Orbitrap Exploris™ 480 mass spectrometer and processed in Proteome Discoverer software using Sequest HT and CHIMERY'S. The results demonstrate a more comprehensive data analysis when using CHIMERY'S: over 80% of all MS2 spectra were matched to one or more peptide precursors and the average number of PSMs per spectrum substantially increases.



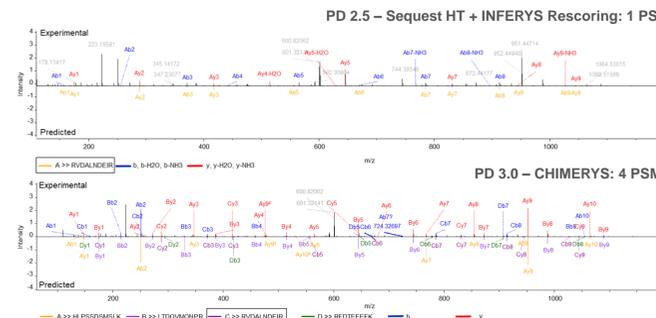
### Figure 3. Number of PSMs per spectrum and identification rate achieved by CHIMERY'S demonstrates the extent of the chimeric spectra problem in DDA data



### Accurate deconvolution by CHIMERY'S unlocks information hidden in chimeric spectra

CHIMERY'S deconvolutes MS2 spectra by considering all peptides for a given spectrum simultaneously, which then compete for the observed experimental intensity in a single step. This results in the identification of several PSMs from chimeric spectra. Using the Proteome Discoverer Spectrum Viewer functionality with direct connection to INFERY'S 2.0, users can visualize the proportional contributions of the individual peptides for every single MS2 spectrum in a mirror plot.

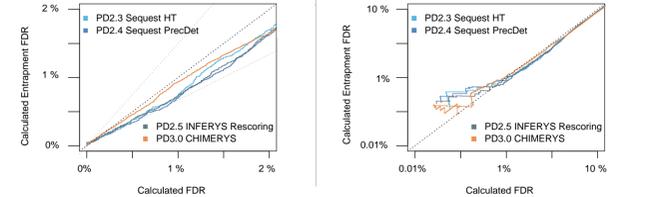
### Figure 5. Mirror plot of an experimental spectrum and PSMs identified by Sequest HT and INFERY'S Rescoring (top panel) or CHIMERY'S (bottom panel) at 1% FDR. While INFERY'S Rescoring identifies only one peptide, CHIMERY'S identifies three additional peptides, resulting in a drastically increased explained intensity of the experimental spectrum.



### Validation of CHIMERY'S results using entrapment searches

Double-decoy approaches enable the calculation of an entrapment FDR and are common benchmarking methods to determine the correctness of FDR estimations. Here, we utilized a human database and appended 8 different plant databases (ratio ~1:7.5 proteins; shared peptides including I/L isomers were removed) to demonstrate the accuracy of the PSM-level FDR calculation performed by Percolator on CHIMERY'S search results of a 1h HeLa cell lysate measurement.

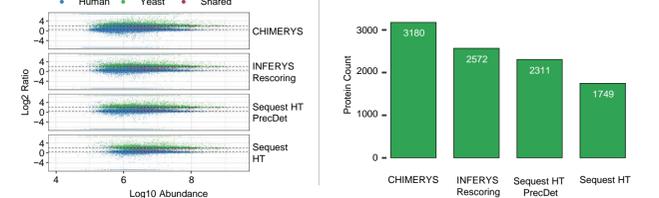
### Figure 6. Analysis of entrapment FDR and calculated FDR using a ~8x non-homologous plant database across different search engines and workflows



### CHIMERY'S increases the number of accurately quantifiable peptides

Due to the increased analysis depth and comprehensive identification of PSMs and peptides, CHIMERY'S aids in the accurate quantification of label-free datasets. We demonstrate this using a two-organism dilution series and compare the quantification results using the Minora feature detector node. This demonstrates that CHIMERY'S produces more quantified peptides and proteins, especially low abundant ones. In this case, CHIMERY'S quantifies 1.8-fold more proteins compared to Sequest HT.

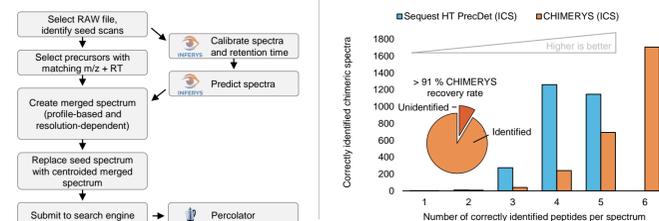
### Figure 8. Quantification of peptide ratios from a HeLa/Yeast dilution experiment (125ng/250ng)



### CHIMERY'S demonstrates an exquisite sensitivity in simulation experiments

To validate CHIMERY'S, we developed an *in-silico* chimeric spectra system (ICS) that spikes *in-silico* generated chimeric spectra into raw files, which can then be used as a ground truth dataset to evaluate search algorithms. Briefly, the system selects seed MS2 spectra with high-confidence identifications from a prior database search from a raw file and convolutes them with several predicted MS2 spectra. To create realistic chimeric data, predicted spectra are derived from peptides with a precursor m/z value within the isolation window of the seed MS2 spectrum and a similar predicted retention time. The created raw file is then submitted to both CHIMERY'S and Sequest HT. Using this system, we demonstrate the sensitivity of CHIMERY'S, which recovers >91% of the *in-silico* chimeric spectra in the convoluted data.

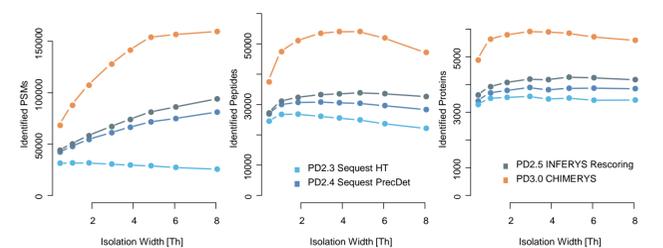
### Figure 10. Schema of the ICS system for generating a ground-truth dataset containing *in-silico* chimeric spectra



### CHIMERY'S enables optimized acquisition settings and profits from increased MS2 complexity

CHIMERY'S deconvolution algorithm is compatible with highly complex samples resulting in convoluted MS2 spectra. Hence, it allows for optimizing data acquisition settings to increase measurement efficiency by identifying more proteins per unit time. Here, we demonstrate that CHIMERY'S enables wider DDA isolation windows that result in more chimeric MS2 spectra, providing more identifications and better sequence coverage.

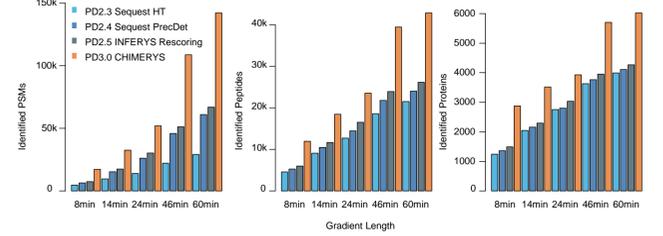
### Figure 12. Number of PSMs, peptide and protein groups identified from a DDA HeLa cell lysate digest acquired on a Thermo Scientific™ Orbitrap Eclipse™ Tribrid™ mass spectrometer using a 1-hour gradient and MS2 isolation windows between 0.4 Th and 8 Th.



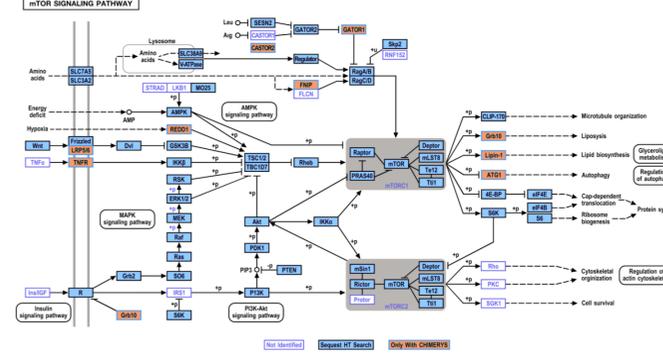
### CHIMERY'S enables increasing sample throughput by using shorter chromatography

CHIMERY'S uniquely deciphers complex samples and MS2 spectra, enabling shorter gradients for LC-MS/MS measurements without losing peptide or protein information in comparison to Sequest HT. Here, we demonstrate how CHIMERY'S identifies the same number of peptides and protein groups in 1/3 of the measurement time.

### Figure 13. Number of PSMs, peptide and protein groups identified by CHIMERY'S or Sequest HT from digests of a HeLa cell lysate acquired on an Orbitrap Exploris 480 MS with gradient lengths ranging from 8 to 60 minutes on a Thermo Scientific™ Vanquish™ Neo UHPLC system



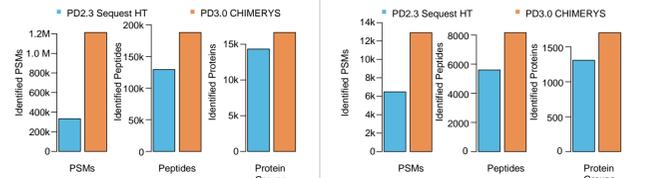
### Figure 14. Proteins of the mTOR signaling pathway identified by CHIMERY'S or Sequest HT in a HeLa cell lysate demonstrate the potential for new biological insight through extended protein and pathway coverage that can be generated using CHIMERY'S



### CHIMERY'S outperforms Sequest HT on datasets from different biological sources

CHIMERY'S is fueled by predictions from INFERY'S 2.0 that are independent of the sample source under investigation. Paired with its resilience with respect to highly complex data, CHIMERY'S is well-equipped to handle fractionated or non-fractionated measurements from organisms from all kingdoms of life [4] and less complex samples like body fluids [5]. Here, we demonstrate its capabilities on a selection of publicly available data.

### Figure 15. Protein groups identified by CHIMERY'S and Sequest HT for a fractionated Arabidopsis thaliana proteome; raw data from PRIDE Project PXD019483 [4]



## CONCLUSIONS

- CHIMERY'S is an innovative, cloud-native search algorithm that uses AI-based predictions to deconvolute chimeric spectra and is fully integrated into Proteome Discoverer 3.0 software
- Using CHIMERY'S results in drastically increased numbers of PSM, peptide and protein group identifications, higher sequence coverage and more confident quantification
- CHIMERY'S excels at analyzing complex samples, enabling more efficient measurements, advanced acquisition settings and shorter gradients to enhance proteomic throughput, productivity and efficiency

## REFERENCES

- Zolg, DP; Gessulat, S; Paschke, C; Frejno M, et al. INFERY'S rescoring: Boosting peptide identifications and scoring confidence of database search results. *Rapid Commun Mass Spectrom*. 2021;e9128. <https://doi.org/10.1002/rcm.9128>
- Dorfer V; Maltsev S; Winkler S; Metchler K. CharmerT: Boosting Peptide Identifications by Chimeric Spectra Identification and Retention Time Prediction. *Journal of Proteome Research* 2018 17 (8), 2581-2589. <https://doi.org/10.1021/acs.jproteome.7b00836>
- The M; MacCoss MJ; Noble WS; Käll L. Fast and Accurate Protein False Discovery Rates on Large-Scale Proteomics Data Sets with Percolator 3.0. *J Am Soc Mass Spectrom*. 2016;27(11):1719-1727. <https://doi.org/10.1007/s13361-016-1460-7>
- Müller JB; Geyer, PE; Colaco, A.R; Mann M; et al. The proteome landscape of the kingdoms of life. *Nature* 582, 592-596 (2020). <https://doi.org/10.1038/s41586-020-2402-x>
- Bian, Y; Zheng, R; Bayer, FP; Kuster B; et al. Robust, reproducible and quantitative analysis of thousands of proteomes by micro-flow LC-MS/MS. *Nat Commun* 11, 157 (2020). <https://doi.org/10.1038/s41467-019-13973-x>

## ACKNOWLEDGEMENTS

The authors would like to thank the alpha and beta testers and especially Karl Metchler for the evaluation and validation of CHIMERY'S. The authors wish to thank numerous scientists and colleagues at MSAID and Thermo Fisher Scientific for fruitful discussions and technical assistance.

## TRADEMARKS/LICENSES

© 2021 Thermo Fisher Scientific Inc. All rights reserved. CHIMERY'S, INFERY'S and MSAID are trademarks of MSAID GmbH. SEQUEST is a trademark of the University of Washington. Powered by AWS is a trademark of Amazon, Inc. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.